

PRODUCT INFORMATION



2''-O-Rhamnosylcariside II

Item No. 36487

CAS Registry No.: 135293-13-9
Formal Name: 3-[[6-deoxy-2-O-(6-deoxy- α -L-mannopyranosyl)- α -L-mannopyranosyl]oxy]-5,7-dihydroxy-2-(4-methoxyphenyl)-8-(3-methyl-2-buten-1-yl)-4H-1-benzopyran-4-one

Synonym: 2''-O-Rhys II

MF: C₃₃H₄₀O₁₄

FW: 660.7

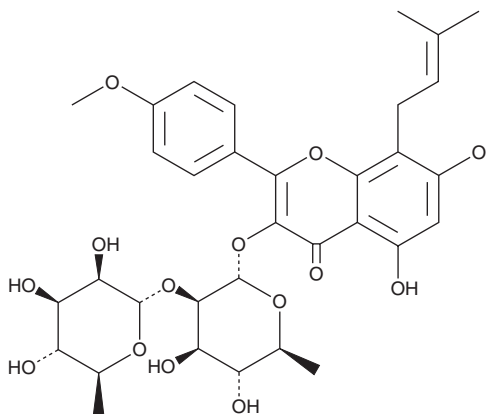
Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Epimedium brevicornu*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

2''-O-Rhamnosylcariside II (2''-O-Rhys II) is supplied as a solid. A stock solution may be made by dissolving the 2''-O-Rhys II in the solvent of choice, which should be purged with an inert gas. 2''-O-Rhys II is soluble in DMSO.

Description

2''-O-Rhys II is a flavonoid glycoside that has been found in *Epimedium* and has hepatotoxic activity.^{1,2} It increases aspartate aminotransferase (AST), lactate dehydrogenase (LDH), and malondialdehyde (MDA) levels, as well as decreases glutathione (GSH) levels in HL 7702 cells when used at a concentration of 66 $\mu\text{g}/\text{ml}$. 2''-O-Rhys II (100 μM) induces 100% mortality in zebrafish larvae.² It also induces hepatocellular vacuolization and degeneration in adult zebrafish when used at a concentration of 25 μM .

References

1. Zhang, L., Wang, T., Zhao, B.-S., et al. Effect of 2''-O-rhamnosyl icaraside II, baohuoside I and baohuoside II in *Herba epimedii* on cytotoxicity indices in HL-7702 and HepG2 cells. *Molecules* **24**(7), 1236 (2019).
2. Zhong, R., Chen, Y., Ling, J., et al. The toxicity and metabolism properties of *Herba epimedii* flavonoids on larval and adult zebrafish. *Evid. Based Complement. Alternat. Med.* 3745051 (2019).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/24/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM